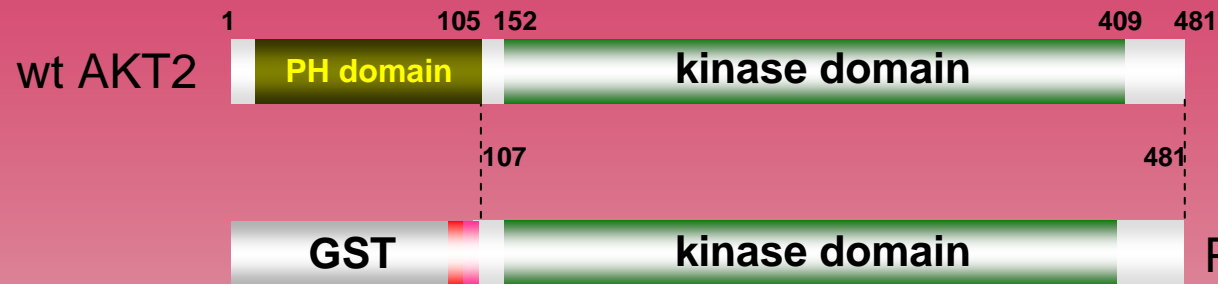


human **AKT2** kinase (GenBank Acc.No: NM_001626.2)



ProQinase AKT2 (Ala₁₀₇-Glu₄₈₁)

633 aminoacids, 72.844 kDa

1-218

GST

224-229

HIS-tag

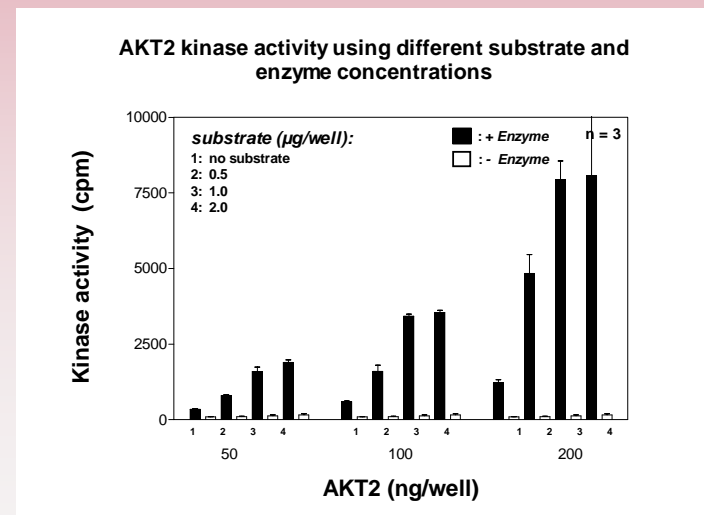
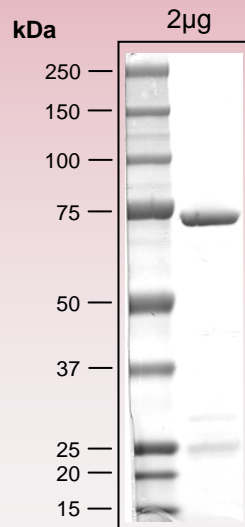
241-246

Thrombin cleavage site

259-633

AKT2 107-481

Representative lot: Coomassie stain & activity in PanQinase™ assay



Lot specific datasheet
available on request.
please contact:
info@proqinase.com